

Data Sheet

MAOA

Human, recombinant, N-terminal His-tag,
N-terminal FLAG-tag

Catalog #: 71103

Lot #: 130822-HF **Conc.:** 0.05 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, 80 µg/ml FLAG peptide, and 20%
glycerol.

Stability: At least 6 months at -70°C. *Avoid
freeze/thaw cycles.*

References:

1. Hodorova, I., *et al.*, *Med. Sci. Monit.*
2012;**18(12)**:BR482-BR486
2. Brunner, H.G., *et al.*, *Science.*
1993;**262(5133)**:578-580.

Description:

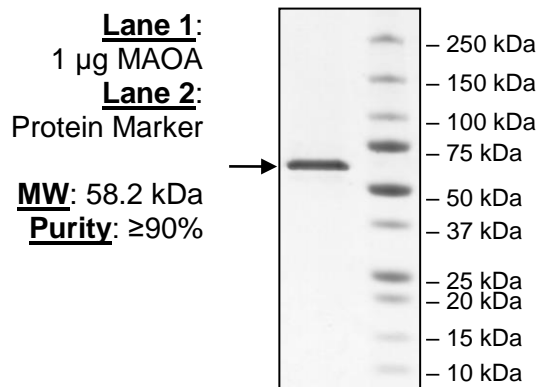
Human monoamine oxidase A (MAOA),
GenBank Accession No. NM_000240, a.a.
2-497, with N-terminal His-tag and N-
terminal FLAG-tag, MW = 58.2 kDa,
expressed in a baculovirus infected Sf9
cell expression system.

Application:

Suitable for protein binding studies, as an
immunogen, or as a protein standard for
Western blotting.

Quality Assurance

4-20% SDS-PAGE Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MHHHHHHDYKDDDDKENQEKASIAGHMFVDDVIGGGISGLSAAKLLTEYGVSVLVLEA
RDRVGGRTYTIRNEHVDYVDVGGAYVGPTQNRILRLSKELGIETYKVNVSERLVQYVKG
KTYPRGAFPPVWNPIAYLDYNNLWRTIDNMGKEIPTDAPWEAQHADKWDKMTMKELI
DKICWTKTARRFAYLNVNINVTSEPHEVSALWFLWYVKQCGGTTTRIFSVTNGGQERKFV
GGSGQVSERIMDLLGDQVKNHPVTHVDQSSDNIIETLNHEHYECKYVINAIPPTLTAKI
HFRPELPAERNQLIQRLPMGAVIKCMMYYKEAFWKKKDYCGCMIEDEDAPISITLDDTK
PDGSLPAIMGFILARKADRLAKLHKEIRKKKICELYAKVLGSQEALHPVHYEKNWCEE
QYSGGCYTAYFPPGIMTQYGRVIRQPVGRIFAGTETATKWSGYMEGAVEAGERAARE
VLNGLGKVTEKDIWVQEPESKDVP AVEITHTFWERNLPS

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com